



Optimized analysis for yeast and other small cells.

Cellometer® X2 fluorescent cell viability counter

The affordable option for yeast viability, vitality and cell concentration assays for breweries.

- 5 imaging modes to provide multiple assays for quality control of brewing yeasts
- Fluorescence yeast viability and cell concentration assay
- Fluorescence yeast vitality assay
- Yeast cell cycle analysis
- De-cluster software for clumpy yeast
- External computer for more convenient sample ID input, data management
- Single fluorescence model, Cellometer X1, also available



Feature	Cellometer X1	Cellometer X2
Bright field imaging mode	X	X
Single fluorescence imaging mode	X	X
Dual fluorescence imaging mode		X
Viability using PI	X	X
Concentration & viability using AOPI		X
Vitality using CFDA-AM		X
Yeast cell cycle	X	X
Cellometer software for analysis of clumpy and irregular-shaped cells	X	X
Mean diameter and cell size distribution	X	X
Cell type wizard for creating new cell type parameters	X	X

For research use only. Not approved for diagnostic or therapeutic use.